# **PPACK**

#### Thrombin inhibitor

Potent irreversible thrombin inhibitor. Inhibits thrombin activation of progelatinase A. Pretreatment with PPACK inhibits thrombin-induced cerebral thromboembolism *in vivo*.

Citations: 7

View Online »

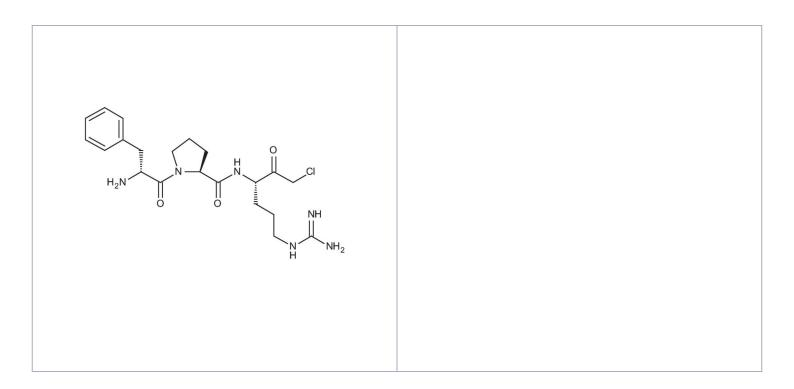
## **Ordering Information**

Order Online »

BML-PI117-0005	5mg
BML-PI117-0025	25mg

Manuals, SDS & CofA

View Online »



## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months

at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name D-Phe-Pro-Arg-CMK

**Appearance** White to off-white solid.

**CAS** 71142-71-7

Formula  $C_{21}H_{31}CIN_6O_3$ 

**MW** 451.0

Purity ≥90% (HPLC)

Sequence D-Phe-Pro-Arg-chloromethylketone

**Solubility** Soluble in water.

**Technical Info / Product Notes**Note: Product is not sterile.

Last modified: May 29, 2024

